

MOUSE RECOMBINANT M4 RECEPTOR MULTISCREEN™ STABLE CELL LINE

Data sheet

PRODUCT INFORMATION

Catalog Number: Cm1025-1

Lot Number: Cm1025-1-100522

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: CellBanker 2

Host cell: CHO-K1

Transfection: Expression vector containing full-length mouse M4 cDNA (GenBank Accession Number NM_007699.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS, 10 ug/mL puromycin

Stability: Stable for a minimum of 2 months in continuous culture

Background: The human muscarinic M4 receptor is a 479-amino acid, 7-transmembrane protein. In heart muscle, M4 receptors are coupled to specific K⁺ currents that regulate cardiac muscle contractions. They can also be found in tissues of the central nervous system and bladder.

Application: Functional assays

Figure 1

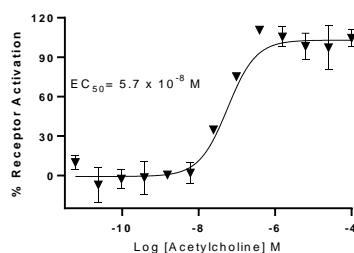


Figure 2

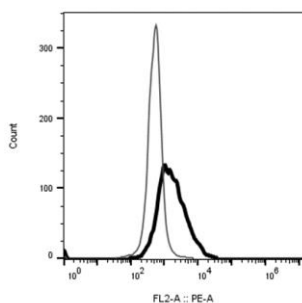


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Eglen (2006) Muscarinic receptor subtypes in neuronal and non-neuronal cholinergic function. *Auton Autacoid Pharmacol* 26:219-233.

Yeh et al. (2007) Atrial tachycardia induces remodeling of muscarinic receptors and their coupled potassium currents in canine left atrial and pulmonary vein cardiomyocytes. *Br J Pharmacol* 152:981-982.

FOR RESEARCH USE ONLY.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.